

## DRI™ TRICYCLICS SERUM TOX APPLICATION



# ORTHO CLINICAL DIAGNOSTICS VITROS® XT 7600 INTEGRATED SYSTEM, VITROS® 5600 INTEGRATED SYSTEM, AND VITROS® 4600 CHEMISTRY SYSTEM

Reference No. 1128

The DRI Tricyclics Serum Tox Assay is intended for the qualitative and semiquantitative determination of tricyclic antidepressants in human serum, plasma, or urine. **This application sheet is for qualitative use only.**



For In Vitro Diagnostic Use Only  
Rx Only

### Intended Use



The information provided in this application sheet is intended as a supplement to the package insert. Refer to the package insert for information on intended use, reagent storage, reagent preparation, specimen collection, specimen preparation, specimen storage, quality control, and additional performance data. For package inserts, visit [www.orthoclinicaldiagnostics.com](http://www.orthoclinicaldiagnostics.com) > Ortho Care > Technical Documents > MicroTip Partnership Assays (MPA).

### Ordering Information

Please place your order with Ortho Clinical Diagnostics. Ordering information available on [www.orthoclinicaldiagnostics.com](http://www.orthoclinicaldiagnostics.com).

### Technical Support Information

Contact Ortho Clinical Diagnostics for technical support. Contact information available on [www.orthoclinicaldiagnostics.com](http://www.orthoclinicaldiagnostics.com).



Microgenics Corporation, part of Thermo Fisher Scientific  
46500 Kato Road, Fremont, CA 94538 USA  
U.S. Toll free: (800) 232-3342 / Tel: (510) 979-5000  
U.S. Toll free fax: (888) 527-8001 / Fax: (510) 979-5420



B.R.A.H.M.S GmbH, Neuendorfstrasse 25, 16761 Hennigsdorf, Germany

*Continued on next page*

## Reagent Pack Storage

Unopened reagents are stable until the expiration date when stored at 2 to 8°C.

Reagents stored in UDxx reagent packs onboard the analyzer are stable for 30 days.

Reagents are supplied liquid ready-to-use.

The UDxx reagent pack may be filled up to 12.25 mL of Reagent A in UDxx/A bottle and 3.75 mL of Reagent B in UDxx/B bottle, which provides 76 tests.

R1 (mL) in UDxx/A	R2 (mL) in UDxx/B	Tests/pack
12.25 mL	3.75 mL	76

**If Splitting:** The provided kit configuration will provide 2 UDxx reagent packs with 152 tests.

**NOTE:** Once the individual UDxx pack number is selected for use during the protocol programming, it is the only UDxx pack number to use for this protocol.

### Special Reagent Packs for User Defined Assays

(Please order from Ortho Clinical Diagnostics; not available from Microgenics)

Part Number	Description	Quantity
680 2246	UD01 Packs (Empty)	1 BOX/6PKS
680 2247	UD02 Packs (Empty)	1 BOX/6PKS
680 2248	UD03 Packs (Empty)	1 BOX/6PKS
680 2249	UD04 Packs (Empty)	1 BOX/6PKS
680 2250	UD05 Packs (Empty)	1 BOX/6PKS
680 2251	UD06 Packs (Empty)	1 BOX/6PKS
680 2252	UD07 Packs (Empty)	1 BOX/6PKS
680 2253	UD08 Packs (Empty)	1 BOX/6PKS
680 2254	UD09 Packs (Empty)	1 BOX/6PKS
680 2255	UD10 Packs (Empty)	1 BOX/6PKS
684 4449	UD11 Packs (Empty)	1 BOX/6PKS
684 4448	UD12 Packs (Empty)	1 BOX/6PKS
684 4445	UD13 Packs (Empty)	1 BOX/6PKS
684 4442	UD14 Packs (Empty)	1 BOX/6PKS
684 4447	UD15 Packs (Empty)	1 BOX/6PKS
684 4444	UD16 Packs (Empty)	1 BOX/6PKS
684 4441	UD17 Packs (Empty)	1 BOX/6PKS
684 4446	UD18 Packs (Empty)	1 BOX/6PKS
684 4443	UD19 Packs (Empty)	1 BOX/6PKS
684 4440	UD20 Packs (Empty)	1 BOX/6PKS
680 2256	UDDL1 Packs (Empty)	1 BOX/6PKS
680 2257	UDDL2 Packs (Empty)	1 BOX/6PKS

---

**Results and  
Data  
Interpretation**

For the Qualitative assay, the VITROS<sup>®</sup> XT 7600 System, VITROS<sup>®</sup> 5600 System, and VITROS<sup>®</sup> 4600 System can report both a unit-less Index number and a NEG / POS classification. The assay is calibrated at the cutoff concentration so that the Index number less than the cutoff value are negative and those greater than the cutoff value are positive.

**NOTE:** *Qualitative results must only be reported as Negative or Positive. Unit-less Index numbers cannot be used in reportable results.*

---

**Calibration  
Frequency**

It is recommended that recalibration occur after a reagent pack change, a calibrator lot change, and as required following quality control procedures.

---

**DRI Tricyclics Serum Tox Assay – Qualitative**  
**Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and**  
**VITROS® 4600 System Parameters**

Full Assay Name: Tricyclics Qual

Version Date: 04/16/2019

Short Assay Name: TCA300

Fluid Type: Serum/Plasma

Assay Model Type: 2 Point Rate

Template: \*2Pt R1-S-R2

Cal Model Type: Linear

Calibrator Bottles: 2

Reagent Reps per Cal: 2

---

**Reagent Lot Information**

**Edit Dilution Parameters**

On-Board Stability: 30 Days

Diluent: DATDIL

Standard Dilution Factor: 1.0

Reagent Lot Num: Kit Lot

Reflex Dilution: OFF

Dilution Factor: 20.0

Shelf Exp. Date: Kit Exp Date

Reduction Factor: 1.0

---

**Edit Result Parameters**

Units: ng/mL

Number of Qualitative Ranges: 2

Significant Digits: 5 Precision Digits: 1

1. NEGATIVE

User Adjusted Parameters

2. POSITIVE Cutoff cal value or greater

Slope: 1.0 Intercept: 0.0

Report Qualitative Result Outside of Range: Yes

CuveTip Exp Time: 35 Temp Sens : No

Measuring (Reportable) 0 to 10000

**(More Assay Params) – Edit 2 Pt Rate Additional Parameters**

Initial Abs. Limits: -0.20 to 3.50

Second Abs. Limits: -0.20 to 3.50

Antigen Excess Factor: 9.0

---

**DRI Tricyclics Serum Tox Assay – Qualitative**  
**Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and**  
**VITROS® 4600 System Parameters, *continued***

**Edit Protocol Parameters**

Step	Volume	Pack ID	Seconds	Wavelength
1. Reagent	126 µL	UDxx /A		
2. Incubation			0.0	
3. Sample	6 µL			
4. Incubation			304.0	
5. Reagent	42 µL	UDxx /B		
6. Incubation			38.0	
7. Read				340 nm
8. Incubation			76.0	
9. Read				340 nm

**Edit Calibration Parameters**

Bottle #	Dil Factor	Cal Rep Resp Range	Calibrator Lot: <u>Cal Kit lot</u>
1	<u>1.0</u>	<u>0.20</u>	Cal value: <u>0</u>
2	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>300</u>

**(More Cal Parm) – Edit Linear or Logit/Log Additional Parameters**

Monotonicity: Increase

Max Resp High: 3.00

Min. Resp. High: 3.00

Cal Fit Goodness Limit: 0.990

Max Resp. Low: -3.00

Min Resp. Low: -3.00

**Edit Triple Read Parameters**

	Reportable Conc.	Triple Read Limit
Reportable Min.:	<u>0</u>	<u>50</u>
Critical Conc.:	<u>300</u>	<u>20.0 %</u>
Reportable Max.:	<u>10000</u>	<u>20.0 %</u>

© 2019 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. VITROS is a registered trademark of Ortho Clinical Diagnostics.

*End*